

### Human IL-7 ELISA Set (without plates) ab47357

画像数 1

#### 製品の概要

製品名	Human IL-7 ELISA Set (without plates)
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 3 pg/ml
検出範囲	6.25 pg/ml - 200 pg/ml
全工程の試験時間	2h 45m
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's IL-7 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-7 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure human IL-7 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>These reagents are intended for research use only, not for diagnostic purposes.</p> <p>The IL-7 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>Recognizes both natural and recombinant human IL-7.</p>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

#### 製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	15 x 96 tests	5 x 96 tests	10 x 96 tests
Chromogen TMB Substrate Solution		6 vials	2 vials	4 vials
Detection biotinylated anti-IL-7 antibody		3 x 0.55ml	1 x 0.55ml	2 x 0.55ml
IL-7 Capture antibody		3 x 0.25ml	1 x 0.25ml	2 x 0.25ml
IL-7 standard	Yellow	15 vials	5 vials	10 vials
Streptavidin-HRP		3 vials	1 vial	2 vials

**関連性** Interleukin 7 (IL-7) is a lymphoid cell growth factor that affects pre-B, pro-B, and early T cells. IL-7 was previously known as pre-B cell growth factor and lymphopoietin 1. IL-7 supports the growth of early B cells from long-term lymphoid bone marrow cultures. It is mitogenic to thymocytes and enhances the response of cells to other stimuli such as polyhydroxyalkanoate (PHA) and concanavalin A (ConA). IL-7 stimulates the proliferation of CD4+/CD8+ cells. The proliferative response of thymocytes to IL-7 is not affected by antibodies to the T cell growth factors such as IL2, IL4 and IL6, suggesting that IL-7 is capable of stimulating T cell proliferation through a pathway independent of the known T cell growth factors. Mature T cells respond to IL-7 and Con A, but not to IL-7 alone. In myeloid lineage cells, IL-7 upregulates the production of pro-inflammatory cytokines and stimulates the tumoricidal activity of monocytes/macrophages. IL-7 is expressed by adherent stromal cells from various tissues.

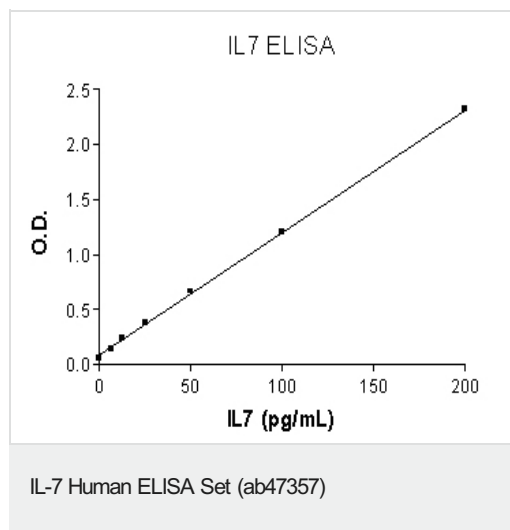
**細胞内局在** Secreted

## アプリケーション

**The Abpromise guarantee** Abpromise保証は、次のテスト済みアプリケーションにおけるab47357の使用に適用されます  
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

## 画像



IL-7 Standard curve ranging from 6.25 to 200 pg/mL

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors